



## Mouse IFN $\gamma$ ELISA

Catalog Number EA-2404

(For Research Use Only)

### Introduction

Interferons (IFNs) are potent extracellular protein mediators of host defence and homeostasis. They are cytokines produced by the cells of the immune system in response to challenges by foreign agents such as viruses, parasites and tumor cells. It is produced by a wide variety of cells in response to the presence of double-stranded RNA, a key indicator of viral infection. IFNs are divided into two major subgroups. Type I IFNs all bind to a type I IFN receptor, such as IFN- $\alpha$  and IFN- $\beta$ . IFN- $\gamma$  is the sole type II IFN, which binds to a distinct type II receptor. Almost all cell types produce type I IFNs, while the type II IFN- $\gamma$  is produced in T cells and natural killer (NK) cells upon immunological stimulation. IFN- $\gamma$  coordinates a diverse array of cellular programs through transcriptional regulation of immunologically relevant genes. Cellular effects of IFN- $\gamma$  includes up-regulation of pathogen recognition, antigen processing and presentation, the antiviral state, inhibition of cellular proliferation and effects on apoptosis, activation of microbicidal effector functions, immunomodulation, and leukocyte trafficking.

### Principle of the assay

IFN $\gamma$  ELISA is based on the principle of a solid phase enzyme-linked immunosorbent assay. The assay utilizes rabbit anti-mouse IFN $\gamma$  antibodies for immobilization on the microtiter wells and rabbit anti-mouse IFN $\gamma$  antibodies along with streptavidin conjugated to horseradish peroxidase (HRP) for detection. The test sample is allowed to react simultaneously with the two antibodies, resulting in the IFN $\gamma$  molecules being sandwiched between the solid phase and enzyme-linked antibodies. After incubation, the wells are washed to remove unbound-labeled antibodies. A HRP substrate, TMB, is added to result in the development of a blue color. The color development is then stopped with the addition of Stop Solution changing the color to yellow. The concentration of IFN $\gamma$  is directly proportional to the color intensity of the test sample. Absorbance is measured spectrophotometrically at 450 nm.

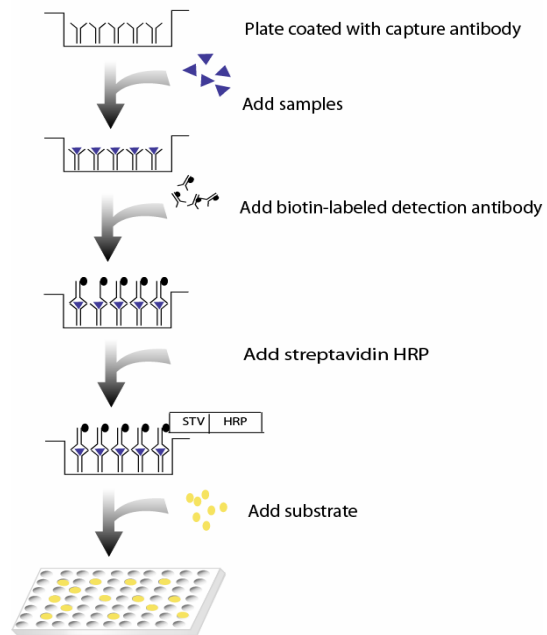


Diagram of ELISA

### Materials provided with the kit

- 96 well microplate coated with rabbit anti-mouse IFN $\gamma$  antibodies (4°C).
- Biotin labeled rabbit anti-mouse IFN $\gamma$  antibodies (-20°C).
- Streptavidin-HRP conjugate (4°C).
- Mouse recombinant IFN $\gamma$  standard (1000ng/ml) (-20°C).
- 1X Diluent buffer (4°C).
- 5X Assay wash buffer (RT)
- Substrate (4°C).
- Stop Solution (4°C).

### Material required but not provided

- Microplate reader capable of measuring absorbance at 450 nm
- Deionized or distilled water.

## Reagent preparation before starting experiment

- Dilute the 5x Assay wash buffer to 1x buffer  
40ml 5x Assay wash buffer  
160ml ddH<sub>2</sub>O
- Dilute 500 times of mouse recombinant IFN $\gamma$  (1000ng/ $\mu$ l) with 1X Diluent blocking buffer to 2000pg/ml and then 2-fold serial dilutions.
- Dilute 400 times of biotin labeled rabbit anti-mouse IFN $\gamma$  antibodies with 1X Diluent buffer before use.
- Dilute 200 times of streptavidin-HRP with 1X Diluent buffer before use.

## Assay procedure

1. Cut the sealing film over the plate and remove it from the desired number of well strips. Make sure the rest of wells are well sealed.
2. Add 100  $\mu$ l of Standard, control, or sample per well and incubate for 1 hour at room temperature with gentle shaking.
3. Aspirate each well and wash by adding 200 $\mu$ l of 1X Assay wash buffer. Repeat the process three times for a total of three washes. Complete removal of liquid at each wash. After the last wash, remove any remaining liquid by inverting the plate against clean paper towels.
4. Add 100 $\mu$ l of diluted biotin-labeled rabbit anti-mouse IFN $\gamma$  antibodies to each well and incubate for 1 hour at room temperature with gentle shaking.
5. Repeat the aspiration/wash as in step 3.
6. Add 100  $\mu$ l of diluted streptavidin-HRP conjugate to each well and incubate for 45 min at room temperature with gentle shaking.
7. Repeat the aspiration/wash as in step 3.
8. Add 100 $\mu$ l substrate to each well and incubate for 5-30 minutes.
9. Add 50 $\mu$ l of Stop solution to each well. The color in the wells should change from blue to yellow.
10. Determine the optical density of each well with a microplate reader at 450 nm within 30 minutes.

## Example of standard curve

